

Product Number:	101160
Lot Number:	501388
Expiration:	12-Dec-2018
Platform(s):	HD-1 & SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent †	101263	81631
Detector Reagent	101264	820604
SBG Reagent	101296	823212
Sample Diluent	103314	818008
Calibrators	101299	80880
Controls	101302 & 101316	80881
SBG Diluent (SR-X)††	103158	N/A ¹
RGP Reagent	103159	N/A ¹

† Follow the Simoa™ Cytokine 3-Plex A Advantage Kit Instructions (KI-0011 - HD1/KI-0101 - SR-X) to prepare the Bead Reagent.

†† Specific for use on the SR-X platform - refer to the Simoa™ Cytokine 3-Plex A SR-X Kit Instructions (KI-0101).

¹ Reagents are not Kit Lot Specific

Release Materials

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80881	0.970	0.733-1.20
Control 2	80881	25.6	18.1-30.9
Panel 1	MF-072115-01	38.9	28.9-56.4
Panel 3	MF-072115-03	1.19	0.901-1.48
Panel 4	MF-072115-04	0.805	0.651-1.12

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80881	0.618	0.426-0.703
Control 2	80881	12.4	9.54-14.5
Panel 1	MF-072115-01	0.898	0.778-1.29
Panel 3	MF-072115-03	1.71	1.46-2.19
Panel 4	MF-072115-04	1.62	1.51-2.26

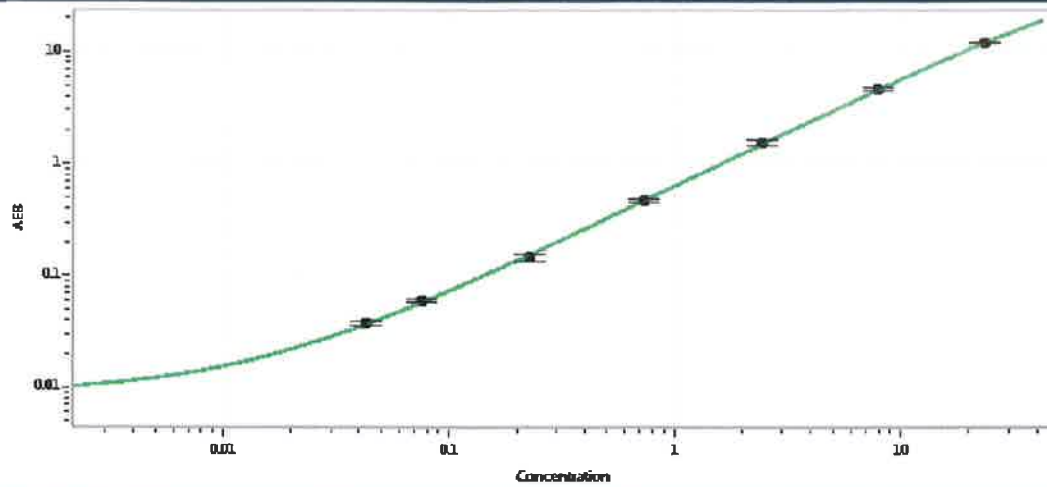
TNFα

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80881	1.90	1.41-2.28
Control 2	80881	47.7	35.1-56.4
Panel 1	MF-072115-01	1.49	0.715-2.32
Panel 3	MF-072115-03	22.4	9.06-34.0
Panel 4	MF-072115-04	6.64	5.43-10.0

Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval	Sivasankari Murugan	QC Supervisor	
	Name	Title	Signature/Date

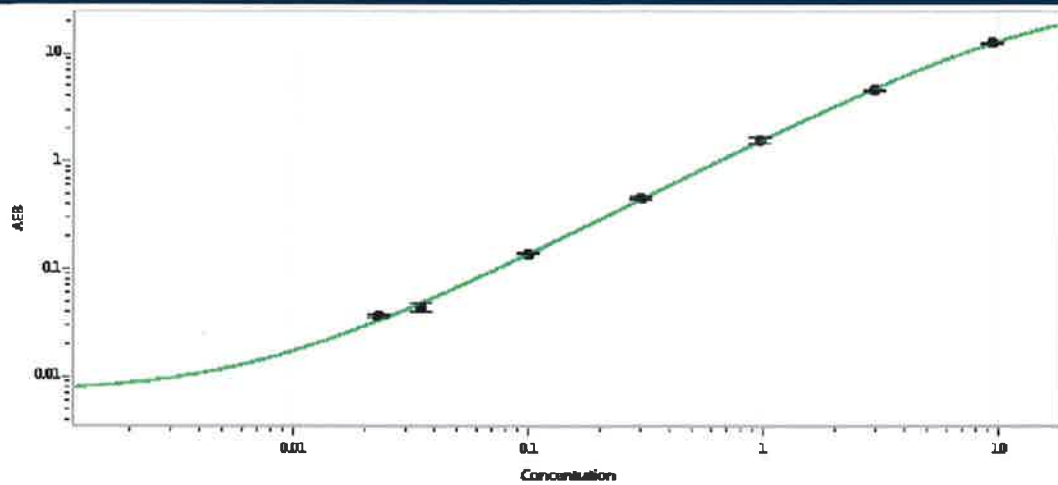
IL-6 Calibration Curve



Calibrator Levels
(pg/mL)

A	0.000
B	0.043
C	0.076
D	0.229
E	0.729
F	2.44
G	7.99
H	23.6

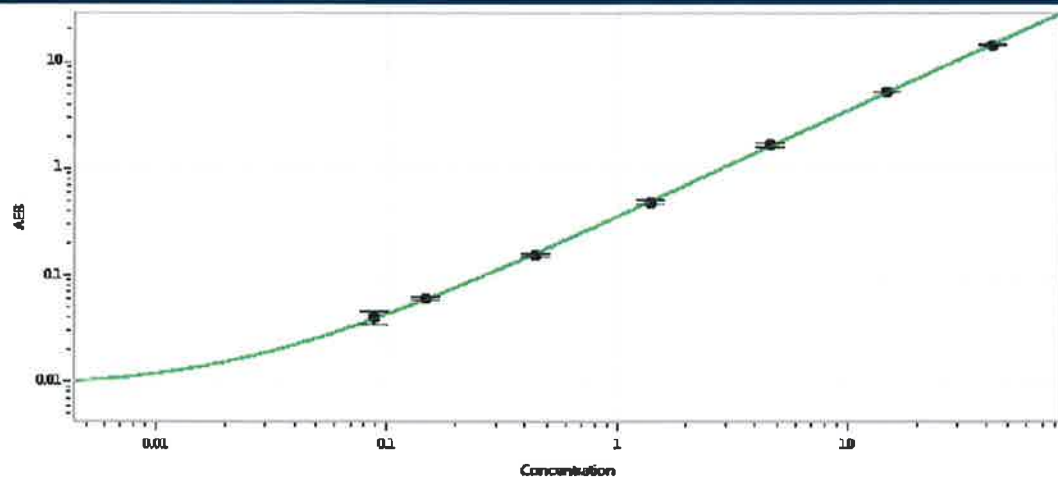
IL-10 Calibration Curve



Calibrator Levels
(pg/mL)

A	0.000
B	0.023
C	0.035
D	0.100
E	0.301
F	0.977
G	2.96
H	9.40

TNFα Calibration Curve



Calibrator Levels
(pg/mL)

A	0.000
B	0.088
C	0.148
D	0.446
E	1.41
F	4.66
G	14.8
H	42.5